

Data Sheet

CD47, Fc-Fusion, Streptavidin-Labeled

Human, Recombinant, Fc fusion protein,
Avi-Tag, Streptavidin-labeled

Catalog #: 71292

Lot#: 170111

Conc.: 3.81 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Vernon-Wilson, E.F. *et al.*, *Eur. J. Immunol.* 2000; **30(8)**: 2130-2137.
2. Majeti, R. *et al.*, *Cell.* 2009; **138(2)**: 286-299.

Description:

Human CD47, Fc fusion protein, also known as leukocyte surface antigen CD47, integrin-associated protein (IAP), OA3, and MER6, GenBank Accession No. NM_001777, a.a. 19-139, with C-terminal Avi-Tag, expressed in a HEK293 cell expression system. The purified protein (40 kDa) was covalently labeled by streptavidin (~13 kDa). Protein runs at a higher MW by SDS-PAGE due to glycosylation and streptavidin labeling.

Application:

Useful as a positive control for western blot analysis.

Quality Assurance

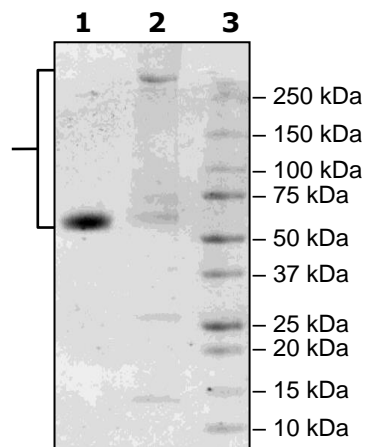
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg CD47 (unlabeled)

Lane 2:
4 μg CD47-Strep

Lane 3:
Protein Marker

MW: CD47 (40 kDa +
glycans) + streptavidin
(monomer = 13 kDa)



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com